
EZH2 Polyclonal Antibody

(Catalog # A-2019)

Background

This gene encodes a member of the Polycomb-group (PcG) family. PcG family members form multimeric protein complexes, which are involved in maintaining the transcriptional repressive state of genes over successive cell generations. This protein associates with the embryonic ectoderm development protein, the VAV1 oncoprotein, and the X-linked nuclear protein. This protein may play a role in the hematopoietic and central nervous systems. Multiple alternatively spliced transcript variants encoding distinct isoforms have been identified for this gene.

Description

EZH2 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.01% thimerosal, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

Q15910

Purification

Affinity Purified

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 50-150 of human EZH2/KMT6 (NP_001190176.1).

Storage

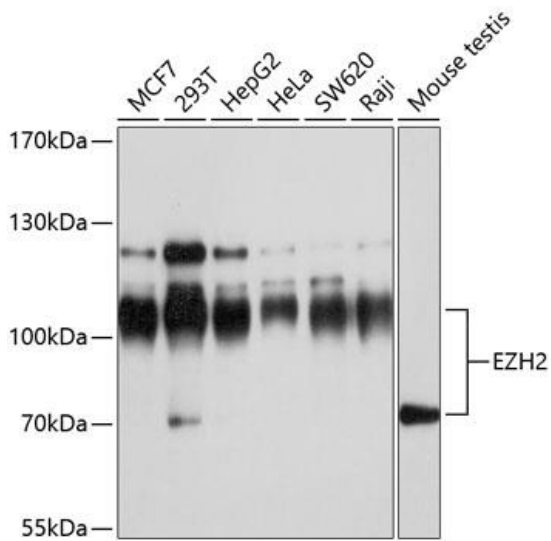
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

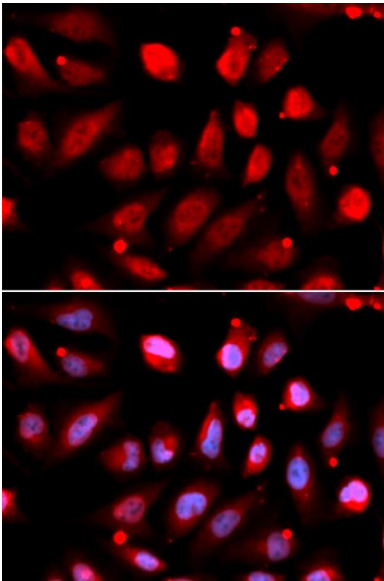
Enhancer of zeste homolog 2 (Drosophila) antibody, ENX 1 antibody, Enx 1h antibody, EZH1 antibody, EZH2b antibody, Histone-lysine N-methyltransferase EZH2 antibody, KMT 6 antibody, KMT6A antibody, Lysine N-methyltransferase 6 antibody, MGC9169 antibody, WVS antibody, WVS2 antibody, WVS; ENX1, EZH1, KMT6, WVS2, ENX-1, EZH2b, KMT6A

Application

WB, IF; Recommended dilution: WB: 1:500-1:2000, IF:1:50-1:200



Western blot analysis of extracts of various cell lines, using EZH2 antibody at 1:1000 dilution.
Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.
Lysates/proteins: 25ug per lane.
Blocking buffer: 3% nonfat dry milk in TBST.



Immunofluorescence analysis of U2OS cells using EZH2 antibody.